

Product Number:	102929
Lot Number:	502950
Expiration:	10-Apr-2022
Platform(s):	HD-1/HD-X



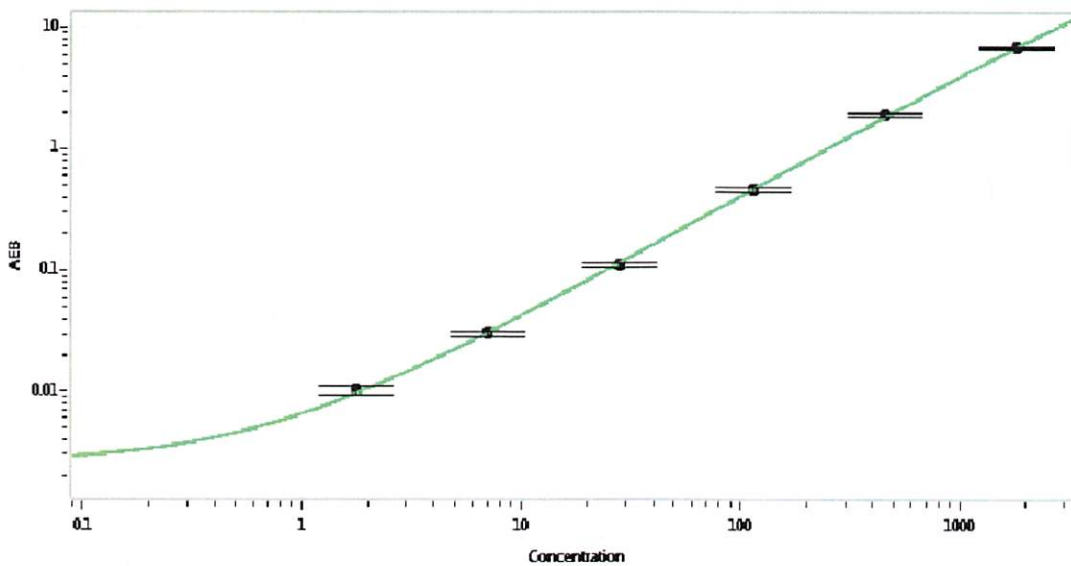
Data below represents results generated on the Simoa™ HD-1 Analyzer.

Component	Part Number	Lot Number
Bead Stock	102922	114109
Detector Stock	102923	116302
Calibrator Concentrate	102926	116303
Calibrator Diluent	102924	114305
Sample Diluent	102925	111023

Generic Components - Included With Kit But Not Lot Number Specific	
Discovery Bead Diluent	101954
Discovery Detector Diluent	101955
Discovery SBG Diluent	101956
Discovery SBG Stock	103297
RGP Reagent	103159

Calibrator Concentrate Target Value	36 ng/mL
-------------------------------------	----------

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	1.76
C	7.03
D	28.1
E	113
F	450
G	1800

Review/Approval

Junpil Park
Name

QC Supervisor
Title

[Signature] 25 Jun 2021
Signature/Date